

DEACM-caged cAMP

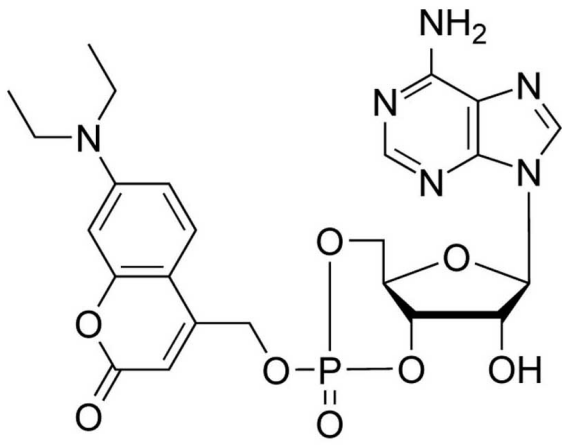
DEACM-caged cAMP is an only weakly fluorescent, caged precursor of the second messenger cyclic AMP, which upon irradiation, releases both cAMP and a fluorescent coumarin analogue. High lipophilicity but still reasonable solubility in aqueous systems.

BLG-D042-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Ordering Information [Order Online »](#)

BLG-D042-001	0.1 μ mol
BLG-D042-005	5x0.1 μ mol

Manuals, SDS & CofA [View Online »](#)



Handling & Storage

Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (7-Diethylaminocoumarin-4-yl)methyl adenosine-3',5'-cyclic monophosphate

CAS 339291-37-1

Couple Type Modified nucleotides

Formula $C_{24}H_{27}N_6O_8P \cdot H_2O$

MW 558.5 . 18.0

Purity >98% by HPLC

Quantity 0.1 μ mol= \sim 58 μ g

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com